

CXXC1 antibody (pAb)

Catalog No: 39203

RRID: AB_2615050

Application(s): WB

Reactivity: Human

Volume: 200 µl

Purification: Protein G Chromatography

Host: Rabbit

Isotype: IgG

Concentration: 0.5 µg/µl

Molecular Weight: 88 kDa

Background: CXXC1 (also known as CGBP or CpG binding protein) is a transcriptional activator protein that binds unmethylated CpG dinucleotides and plays a role in regulating gene expression and embryonic development. CXXC1 exhibits a different distribution pattern than other heterochromatin-associated proteins. It seems to co-localize with acetylated histones, indicating that CXXC1 is associated with active chromatin

Immunogen: This CXXC1 antibody was raised against a mixture of synthetic peptides corresponding to amino acid residues 73-90 and 641-656 of human CXXC1.

Buffer: PBS containing 0.05% BSA and 0.02% sodium azide. Sodium azide is highly toxic.

Application Notes:

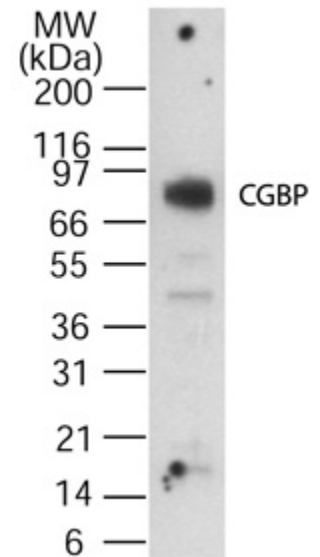
Applications Validated by Active Motif:

WB: 1:500 - 1:1,000 dilution

For optimal results, primary antibody incubations should be performed at room temperature. The addition of 0.1% Tween 20 to all blocking solutions may also reduce background. Individual optimization may be required.

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Store at 4°C for short term. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



CXXC1 pAb tested by Western blot.

Detection of CXXC1 by Western blot analysis using CXXC1 pAb at a 1:1,000 dilution against 15 µg per lane of human spleen lysate.